



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

MTSBS

MTSBS (4-Sulfonatobutyl methanethiosulfonate sodium salt) is a negatively charged MTS reagent.

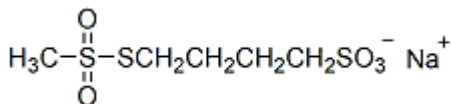


Product attributes

Product Description

MTSBS (4-Sulfonatobutyl methanethiosulfonate sodium salt) is a negatively charged MTS reagent.

- White solid soluble in water, DMSO or DMF
- Store desiccated at -20 °C
- C₅H₁₁NaO₅S₃
- MW: 270



This datasheet was generated on June 22, 2026 at 01:13:55 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mtsbs-4-sulfonatobutyl-methanethiosulfonate-sodium-salt/>